



P.O. Box 58537
Salt Lake City, Ut 84108
Telephone 801-588-0455
Fax 801-588-0497
Email: echelon@echelon-inc.com
http://www.echelon-inc.com

Certificate of Analysis

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Product Name: 1-Deoxy-D-xylulose 5-phosphate (DXP)

Product Number: I-M050

Lot Number: E00162-126

Product Description: Lyophilized powder

Quality Assay: ^1H and ^{31}P NMR

Quality Parameters

A. NMR, Spectra matches control spectra

Passes



Passes All Functional Specifications

A handwritten signature in black ink that reads "Lee Crosby". The signature is written over a horizontal line.

Date 05/06/09

Lee Crosby
Quality Manager